

## Data Sheet

### PRMT8

Human, recombinant, N-terminal His-FLAG-tag  
**Catalog #:** 51052  
 Lot #: 120416                      Conc.: 0.2 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 ng/μl Flag peptide, 3 mM DTT, and 20% Glycerol.

**Stability:** >6 months at -80°C

**References:**

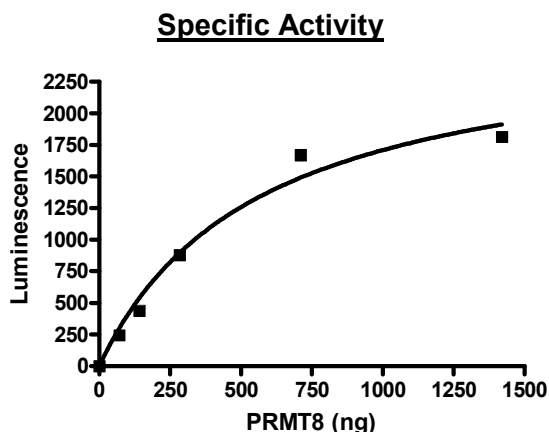
1. Bedford MT., *et al.* J Biol Chem. 2007 Dec 14; **282(50)**:36444-53.
2. Taneda T., *et al.* Brain Res. 2007 Jun 25; **1155**:1-9.

**Description:** Human PRMT8 (protein arginine methyltransferase 8), GenBank Accession No. NM\_019854, a.a. 61-394(end) with N-terminal His-FLAG-tag, MW=40.8 kDa, expressed in Sf9 cells using a Baculovirus expression system.

**Assay Conditions:** 50 μl reaction mix (20 mM phosphate buffer pH7.4, 20 μM S-adenosylmethionine, and 1-5 ng methyltransferase PRMT8) added to the wells coated with the substrate. Incubate at RT for 1 hour. Then, add secondary HRP-labeled antibody and incubate 30 minutes. Finally, add HRP chemiluminescent substrates and read luminescence.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

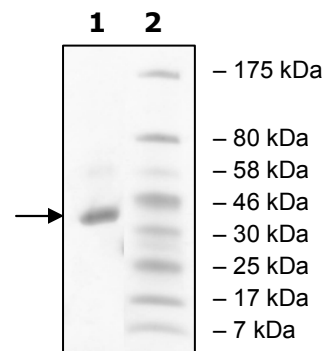
### Quality Assurance



**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2.8 μg PRMT8  
**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 40.8 kDa  
**Purity:** ≥94%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)